

PRODUCT INFORMATION



Benazepril Acyl-β-D-Glucuronide

Item No. 27619

Formal Name: (2S,3S,4S,5R,6S)-2-(2-((S)-3-(((S)-1-ethoxy-1-oxo-4-phenylbutan-2-yl)amino)-2-oxo-2,3,4,5-tetrahydro-1H-benzo[b]azepin-1-yl)acetoxy)-4,5,6-trihydroxytetrahydro-2H-pyran-3-carboxylic acid

Synonym: Benazapril Glucuronide

MF: C₃₀H₃₆N₂O₁₁

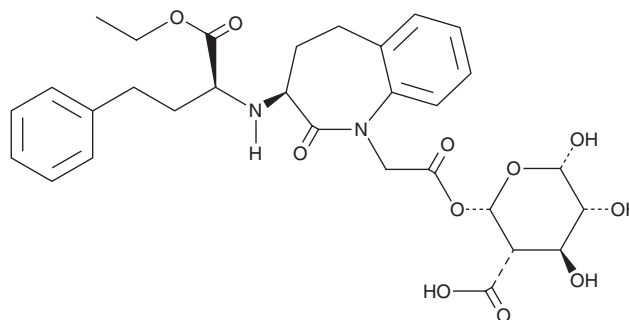
FW: 600.6

Purity: ≥95%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Benazepril acyl-β-D-glucuronide is supplied as a solid. A stock solution may be made by dissolving the benazepril acyl-β-D-glucuronide in the solvent of choice, which should be purged with an inert gas. Benazepril acyl-β-D-glucuronide is slightly soluble in methanol.

Description

Benazepril acyl-β-D-glucuronide is a metabolite of the angiotensin-converting enzyme (ACE) inhibitor prodrug benazepril (Item No. 23804).¹

Reference

1. Xu, L. Benazepril. *Handbook of Metabolic Pathways of Xenobiotics*. Lee, P.W., Aizawa, H., Gan, L.L., Prakash, C., and Zhong, D. editors, 1st edition, John Wiley & Sons, Ltd. (2014).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/19/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM